

Acetylcholine Iodide, [Acetyl- $^3\text{H}$ ]-

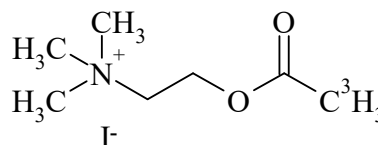
Product Number: NET113

**LOT SPECIFIC INFORMATION**

Lot Number: 3289565

Specific Activity: 83.4 mCi/mmol

Production Date: 20-Jun-2024

M.W. 273.1  
 $\text{C}_7\text{H}_{16}\text{INO}_2$ **PACKAGING:** Crystalline solid in a screw-cap bottle which protects the contents from light.**STABILITY AND STORAGE RECOMMENDATIONS:** When acetylcholine iodide, [acetyl- $^3\text{H}$ ]- is stored at 5°C as a crystalline solid, the rate of decomposition is initially 1% per year from date of purification. The product is not stable in solution; fresh aliquots should be made up immediately prior to use. Stability is nonlinear and not correlated to isotope half-life. Lot to lot variation may occur.**SPECIFIC ACTIVITY RANGE:** 50-100 mCi/mmol (5.46 - 2.73 mg/mCi)**RADIOCHEMICAL PURITY:** This product was initially found to be greater than 97% when determined by the following method. The rate of decomposition can accelerate. It is advisable to check purity prior to use:

Paper chromatography on Whatman No. 1 using the following solvent system:

n-butanol : ethanol : acetic acid : water, (8:2:1:3).

**QUALITY CONTROL:** The radiochemical purity of acetylcholine iodide, [acetyl- $^3\text{H}$ ]- is checked at appropriate intervals using the above chromatography method.**HAZARD INFORMATION:** WARNING: This product contains a chemical known to the state of California to cause cancer.**Limited Warranty:** The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR

FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.

NET113-R Rev02

revvity

Revvity, Inc.  
940 Winter Street  
Waltham, MA 02451 USA

(800) 762-4000  
www.revvity.com

For a complete listing of our global offices, visit [www.revvity.com](http://www.revvity.com)  
Copyright ©2023, Revvity, Inc. All rights reserved.